Biotin-4-fluorescein

Biotin-4-fluorescein (*N*-(5-aminopentyl)biotinamide, trifluoroacetic acid salt) has been reported to be a superior fluorescent biotin derivative for accurate measurement of avidin and streptavidin in crude biological samples. Binding of biotin to avidin and streptavidin is high affinity, fast and noncooperative.

alsitum

Product attributes

 Excitation/Emission
 494/523 nm

 CAS number
 1032732-74-3

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

Biotin-4-fluorescein (*N*-(5-aminopentyl)biotinamide, trifluoroacetic acid salt) has been reported to be a superior fluorescent biotin derivative for accurate measurement of avidin and streptavidin in crude biological samples. Binding of biotin to avidin and streptavidin is high affinity, fast and noncooperative.

- $\lambda_{Ex}/\lambda_{Em} = 494/523 \text{ nm}$
- Yellow solid soluble in DMSO, or pH>7 buffers
- Store at -10 °C to -35 °C and protect from light
- C₃₃H₃₂N₄O₈S
- MW: 644.7

References

- 1. Biochim Biophys Acta 1427, 33 (1999).
- 2. Biochim Biophys Acta 1427, 44 (1999).

This datasheet was generated on November 12, 2025 at 09:15:28 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/biotin-4-fluorescein/